

## Recombinant PRDM5 protein

**Catalog No:** 31494, 31898

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.25 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant PRDM5 protein was expressed in Sf9 cells and is supplied at a concentration of 0.25 µg/µl in 25 mM Hepes pH 7.5, 400 mM NaCl, 5% Glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

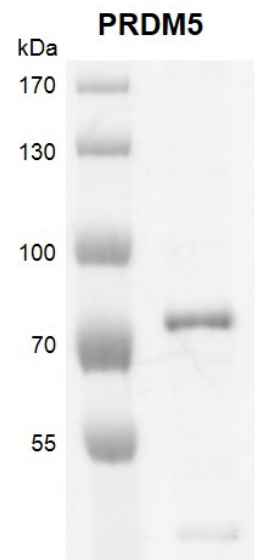
**Background:** **PRDM5** (PR Domain Containing 5) is a transcription factor of the PR-domain protein family. It contains a PR-domain and multiple zinc finger motifs. Mutations of PRDM5 causes Type 2 brittle cornea syndrome (BCS2) which is an inherited connective tissue disease with a devastating ocular phenotype through histone and DNA methylation. PRDM5 is a relevant tumour suppressor gene that is frequently targeted in colorectal tumourigenesis. Promoter methylation-mediated downregulation of PRDM5 contributes to the development of lung squamous cell carcinoma. It is a sequence-specific DNA-binding transcription factor and represses transcription at least in part by recruitment of the histone methyltransferase EHMT2/G9A and histone deacetylases such as HDAC1.

**Protein Details:** Recombinant PRDM5 protein (accession number NP\_061169.2) was expressed in Sf9 cells and contains an N-terminal FLAG tag with a molecular weight of 73.1 kDa. The recombinant protein is >85% pure by SDS-PAGE.

**Application Notes:** Recombinant PRDM5 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**Recombinant PRDM5 protein gel**  
PRDM5 protein was run on an 8% SDS-PAGE gel and stained with Coomassie blue.